

Reference # / No. de référence / Referenz-Nr. _____

Date / Date / Datum _____

Tech / Tech / Techn. _____

Source / Source / Quelle _____

Catalase / Catalase / Katalase : + / - _____

Pigment / Pigment / Pigment _____

Reagent / Réactif / Reagenz /	None / Aucun / Keine										RapID CB Plus Reagent / Réactif RapID CB Plus / RapID CB Plus Reagens					None / Aucun / Keine	Nitrate A & B	H ₂ O ₂	None / Aucun / Keine					
	Yellow, gold, or light yellow-orange / Jaune, or ou légèrement jaune orangé / Gelb, goldgelb oder leichtes gelb-orange				Yellow / Jaune / Gelb				Yellow or yellow-orange / Jaune ou Jaune orangé / Gelb oder gelb-orange			Purple, red, or dark pink / Violacé, rouge ou rose foncé / Dunkelrot, rot oder dunkelrosa					Red or dark red-orange/Rouge ou rouge orangé foncé/Rot oder dunkles rot-orange	Red or Pink / Rouge ou rose / Rot oder rosa	Bubbles / Bulles / Luftblasen	Yellow / Jaune / Gelbes				
Cavity #, No. cavité / Kammer-Nr.	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18						
Test Code / Code du test / Testcode	GLU	SUC	RIB	MAL	aGLU	BGLU	NAG	GLY1	ONPG	PHS	EST	PRO	TRY	PYR	LGLY	LEU	URE	NIT	CAT	PIG				
Value / Valeur / Wert	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2				
Result / Résultat / Ergebnis																								
Value Total / Total des valeurs / Gesamtwert																								

IDENTIFICATION / IDENTIFICATION / IDENTIFIZIERUNG _____



Microcode _____

Riferimento N. / N° de referencia _____

Data / Fecha _____

Tech / Tec _____

Origine / Origen _____

Catalasi / Catalasa: + / - _____

Pigmento / Pigmento _____

Reagente Reactivo	Nessuno / Ninguno										RapID CB Plus Reagent / Reactivo RapID CB Plus					Nessuno/ Ninguno	Nitrate A & B	H ₂ O ₂	Nessuno/ Ninguno																				
	Giallo, oro o giallo-arancio chiaro / Amarillo, dorado o amarillo-naranja claro				Giallo Amarillo						Giallo o giallo- arancio / Amarillo o amarillo- naranja					Porpora, rosso o rosa scuro / Púrpura, rojo o rosa oscuro					Rosso o rosso- arancione scuro / Rojo o rojo- naranja oscuro					Rosso o rosa / Rojo o rosa					Bolle / Burbujas					giallo / amarillo			
Cavità N. / N° de cavidad	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18																					
Codice esame Código de prueba	GLU	SUC	RIB	MAL	aGLU	BGLU	NAG	GLY1	ONPG	PHS	EST	PRO	TRY	PYR	LGLY	LEU	URE	NIT	CAT	PIG																			
Valore / Valor	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2	4	1	2																			
Risultato Resultado																																							
Valore totale Valor total																																							

IDENTIFICAZIONE / IDENTIFICACIÓN



Microcode _____